

Technical Data Sheet

AlphaLISA® SureFire® Ultra™ p-Ribosomal Protein S6 (Ser240/244) Assay Kit

ALSU-PS6R-A500, ALSU-PS6R-A10K, ALSU-PS6R-A50K, ALSU-PS6R-A-HV

Kit-Specificity

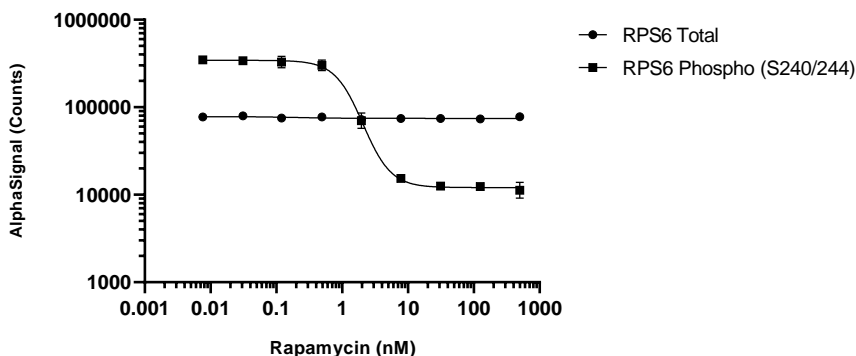
This assay kit contains antibodies which recognize the phospho-RPS6 (Ser240/244) epitope, and a distal epitope, on RPS6. The protein detected by this kit corresponds to UniProt ID P62753. RPS6 is also known as 40S ribosomal protein S6. These antibodies recognize RPS6 of human, mouse and rat origin. Other species should be tested on a case-by-case basis.

Control Lysate Information

Positive Control Lysate: Prepared from MCF7 cells, cultured to confluence in T175 flasks in 10% FBS containing medium, starved for 2 hours, then treated with 500 µg/mL insulin for 60 minutes and lysed in 50 mL of 1X *SureFire Ultra* Lysis Buffer.

Representative Data

Data obtained with a 2-plate, 2-incubation protocol. MCF7 cells were plated at 40K cells per well in a 96 well plate and incubated overnight. Cells were treated with rapamycin at the indicated concentrations for 2 hours prior to treatment with 2.5 µg/mL insulin for 30 minutes. Cells were lysed with 1x Lysis Buffer and assayed for Total or Phospho (Ser240/244) RPS6 using respective *SureFire Ultra* kits. Equivalent to approximately 2000 cells/datapoint.



For more information and detailed protocols, please refer to the generic AlphaLISA *SureFire Ultra* Assay Kits Manual, available from <http://www.perkinelmer.com/category/alpha-surefire-kits>